



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Optimizer Buffer™ -VII

(Cat. #BKC-10)



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Optimizer Buffers™ provide the optimal conditions for specific protein reactions, including protein labeling, modification and cross-linking reactions.

Optimizer Buffer™-VII is designed to carry out maleimide conjugation reaction.

## ITEM(S) SUPPLIED (Cat. # BKC-10)

| Description           | Size      |
|-----------------------|-----------|
| Optimizer Buffer™-VII | 2 x 125ml |

## STORAGE CONDITIONS

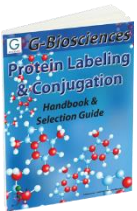
It is shipped at ambient temperature. Store refrigerated at 4°C. This product is stable for 1 year.

## INSTRUCTIONS FOR USE

Optimizer Buffer™-VII is in ready to use form. Simply dissolve sulfhydryl-peptide for maleimide conjugation in Optimizer Buffer™-VII.

## RELATED PRODUCTS

Download our Protein Labeling & Conjugation Handbook.



<http://info2.gbiosciences.com/complete-protein-labeling-conjugation-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

*This page is intentionally left blank.*



[www.GBiosciences.com](http://www.GBiosciences.com)